

Product Number:	101362
Lot Number:	234311
Expiration:	9-Dec-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Khilka

QC Manager

Sonia 12/16/21 DEC 2022

Name

Title

Signature/Date